

Product Number:	102040
Lot Number:	503790
Expiration:	8-Sep-2023

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	234016
Biotinylation Reaction Buffer	101358	234206
Bead Conjugation Buffer	101357	233907
Bead Blocking Buffer	101356	227910
Bead Wash Buffer	101355	225111
Bead Diluent Buffer	100458	232109

Review/Approval

Sonia Khillan

Name

### QC Manager

Title

50112012 / 25 JAN 2023

Signature/Date